

Tecan Award 2011 winner announced



Established in 2010 to celebrate the innovation and ingenuity of Tecan's loyal customers, the Tecan Award highlights some impressive and imaginative uses for the Company's detection instruments. We are pleased to announce that the 2011 award has been won by Svenja Kristina Holle from the University Hospital of Muenster (UHM), Germany, for her work with an Infinite® 200 microplate reader. Svenja and the UHM team have developed an *ex vivo* method for investigating the dynamics of organic cation transport in proximal tubules of the kidney. This innovative optical method could be adapted to almost any kind of freshly isolated biological transport system using a suitable fluorescent substrate, offering an exciting new technique to study membrane transport.

Second place in the 2011 competition went to Michael Bunge from the Institute of Applied Microbiology at Giessen University, Germany, who is using an Infinite M200 to monitor the effects of a metallic nanoparticle-based approach to eliminating the growth of potential pathogenic microorganisms in drinking water (featured on page 30 of this issue). And a close third place went to María Isabel Pividori from the Universitat Autònoma de Barcelona (UAB), Spain, for developing a novel method for the quantification of gliadin in gluten-free foods using a Sunrise™ reader (see page 20 for more information).

The 2011 award will be presented to Svenja during a tour of Tecan's Detection Headquarters in Grödig, Austria, as part of her prize of a trip to the romantic city of Salzburg.

To find out more about the Tecan Award and details of the winning entries, visit www.tecan.com/award



Second Place: Michael Bunge (third from left) from the Institute of Applied Microbiology at Giessen University, Germany, see feature page 30



Third Place: María Isabel Pividori (far right) from the Universitat Autònoma de Barcelona (UAB), Spain, see feature page 20